

## ProQinase™ PIK3CB E633K/PIK3R1

phosphoinositide-3-kinase, catalytic, beta polypeptide/  
phosphoinositide-3-kinase, regulatory subunit 1 (alpha)

### Recombinant Human Active Lipid Kinase

**HGNC Symbol:** PIK3CB

**Synonyms PIK3CB:** P110BETA, PI3K, PI3KBETA, PI3K-beta, PIK3C1

**Synonyms PIK3R1:** GRB1, p85, p85-ALPHA, PtdIns-3-kinase regulatory subunit p85-alpha

### Lipid Kinase Family: PI3K Class I

(according to: Phylogenomics of phosphoinositide lipid kinases: perspectives on the evolution of second messenger signaling and drug discovery: James R Brown & Kurt R Auger; BMC Evolutionary Biology 11, 4-14 (2011))

**Product No.:** 1531-1165-1

**Lot:** 003

**Description:** Human PIK3CB, full length, amino acids M<sub>1</sub>-S<sub>1070</sub> (as in NCBI/Protein entry NP\_006210.1) with a E633K mutation, N-terminal GST-HIS<sub>6</sub> fusion protein with a 3C cleavage site and PIK3R1 full length, amino acids M<sub>1</sub>-R<sub>724</sub> (as in NCBI/Protein entry NP\_852664.1), N-terminal fused to a MYC-tag, expressed in Sf9 insect cells

**Product identity:** PIK3CB E633K/PIK3R1 Lot 003, was confirmed as PIK3CB/PIK3R1 by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW<sub>GST-PIK3CB E633K</sub>:** 151,266 Da

**Theoretical MW<sub>PIK3R1</sub>:** 85,371 Da

**Expression:** Baculovirus infected Sf9 cells

**Purification:** GST-Affinity Chromatography

**Storage buffer:** 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 20% glycerol

**Storage temperature:** -80°C

Avoid repeated freeze-thaw cycles!

For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

**Protein concentration:** 0.116 µg/µl

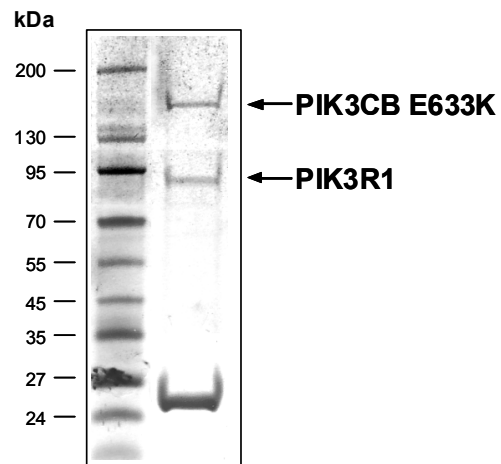
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

### Biochemical Parameters:

Specific kinase activity (P<sub>i</sub> transfer): 78 pmol/µg×min

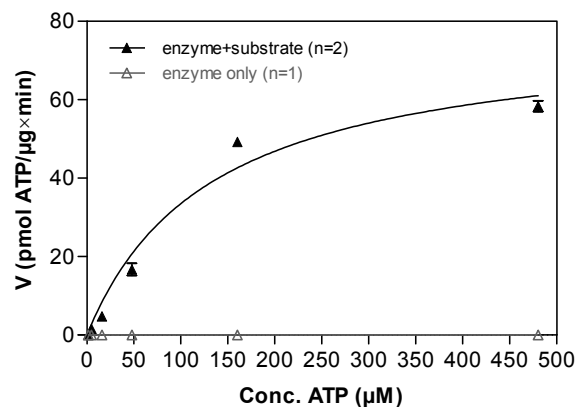
ATP-K<sub>M</sub>: 132 µM

### PIK3CB E633K/PIK3R1 Lot 003: Coomassie stain



2 µg PIK3CB E633K/PIK3R1

### PIK3CB E633K/PIK3R1 Lot 003: Determination of V<sub>max</sub> and K<sub>M</sub> value for ATP ADP-Glo™ Kinase Assay / Promega



### Determination of K<sub>M</sub> value & Specific activity:

• Assay conditions:

60 mM HEPES-NaOH, pH 7.5

3 mM MnCl<sub>2</sub>

3 µM Na-orthovanadate

1.2 mM DTT

50 µg / ml PEG<sub>20,000</sub>

ATP (variable)

Substrate: PIP2: 50 µM / PS: 950 µM

PIP2: 08:0 PI(4,5)P2 (1,2-Dioctanoyl-sn-Glycero-3-(Phosphoinositol-4,5-Bisphosphate))

PS: 1-Palmitoyl-2-Oleoyl-sn-Glycero-3-[Phospho-L-Serine])

PIK3CB E633K/PIK3R1: 4.0 µg / ml

For further information on ADP-Glo™ kinase activity detection please visit [Promega.com](http://www.promega.com)

# ProQinase™ PIK3CB E633K/PIK3R1

Product No.: 1531-1165-1

PIK3CB E633K Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMG <b>HHHHHG</b>	RDS <b>LEVLFOG</b>	240
241	<b>PLAMVMCFSF</b>	<b>IMPPAMADIL</b>	<b>DIWAVDSQIA</b>	<b>SDGSIPVDFL</b>	<b>LPTGIYIQLE</b>	<b>VPREATISYI</b>	300
301	<b>KQMLWKQVHN</b>	<b>YPMFNLLMDI</b>	<b>DSYMFACVNO</b>	<b>TAVYEELEDE</b>	<b>TRRLCDVRPF</b>	<b>LPVLKLVTRS</b>	360
361	<b>CDPGEKLDISK</b>	<b>IGVLIGKGLH</b>	<b>EFDLSKDPEV</b>	<b>NEFRRKMRKF</b>	<b>SEEKILSLVG</b>	<b>LSWMDWLKQT</b>	420
421	<b>YPPEHEPSIP</b>	<b>ENLEDKLYGG</b>	<b>KLIVAVHFEN</b>	<b>CQDVFSFQVS</b>	<b>PNMNPIKVNE</b>	<b>LAIQKRLTIH</b>	480
481	<b>GKEDEVSPYD</b>	<b>YVLQVSGRVE</b>	<b>YVFGDHPLIQ</b>	<b>FQYIRNCVMN</b>	<b>RALPHFILVE</b>	<b>CKKIKKMYEQ</b>	540
541	<b>EMIAIEAAIN</b>	<b>RNSSNLPLPL</b>	<b>PPKKTRIISH</b>	<b>VWENNNPFQI</b>	<b>VLVKGKLNLT</b>	<b>EETVKVHVRA</b>	600
601	<b>GLFHGTELLC</b>	<b>KTIVSSEVSG</b>	<b>KNDHIWNEPL</b>	<b>EFDINICDLP</b>	<b>RMARLCFAVY</b>	<b>AVLDKVKTKK</b>	660
661	<b>STKTINPSKY</b>	<b>QTIRKAGKVH</b>	<b>YPVAWVNTMV</b>	<b>FDKFGQLRTG</b>	<b>DIILHSWSSF</b>	<b>PDELEEMLNP</b>	720
721	<b>MGTVQTNPYT</b>	<b>ENATALHVKF</b>	<b>PENKKQPYYY</b>	<b>PPFDKIIEKA</b>	<b>AEIASSDSAN</b>	<b>VSSRGGKKFL</b>	780
781	<b>PVLKEILDRLD</b>	<b>PLSQLCENEM</b>	<b>DLIWTLRQDC</b>	<b>REIFPQSLPK</b>	<b>LLLSIKWNKL</b>	<b>EDVAQLQALL</b>	840
841	<b>QIWPKLPPRE</b>	<b>ALELLDFNYP</b>	<b>DQYVREYAVG</b>	<b>CLRQMSDKEL</b>	<b>SOYLLQLVQV</b>	<b>LKYEPFLDCA</b>	900
901	<b>LSRFLLERAL</b>	<b>GNNRIGQFLF</b>	<b>WHLRSEVHIP</b>	<b>AVSVQFGVIL</b>	<b>EAYCRGSVGH</b>	<b>MKVLKQVEA</b>	960
961	<b>LNLKLTLSNL</b>	<b>IKLNAVKLNLR</b>	<b>AKGKEAMHTC</b>	<b>LKQSAYREAL</b>	<b>SDLQSPLNPC</b>	<b>VILSELYVEK</b>	1020
1021	<b>CKYMDSKMKP</b>	<b>LWLVIYNNKVF</b>	<b>GEDSVGVIFK</b>	<b>NGDDLQDML</b>	<b>TLQMLRLMDL</b>	<b>LWKEAGLDR</b>	1080
1081	<b>MLPYGCLATG</b>	<b>DRSGLIEVVS</b>	<b>TSETIADIQL</b>	<b>NSSNVAAAAA</b>	<b>FNKDALLNWL</b>	<b>KEYNSGDDLD</b>	1140
1141	<b>RAIEEFTLSC</b>	<b>AGYCVASYVL</b>	<b>GIGDRHSDNI</b>	<b>MVKKTGQLFH</b>	<b>IDFGHILGNF</b>	<b>KSKFGIKRER</b>	1200
1201	<b>VPFILTYDFI</b>	<b>HVIQQGKTGN</b>	<b>TEKFGFRFQC</b>	<b>CEDAYLILRR</b>	<b>HGNLFITLFA</b>	<b>LMLTAGLPEL</b>	1260
1261	<b>TSVKDIQYLK</b>	<b>DSLALGKSEE</b>	<b>EALKQFKQKF</b>	<b>DEALRESWTT</b>	<b>KVNWMAHTVR</b>	<b>KDYRS</b>	1320

1-218: GST    **Red:** HIS6-tag    **Green:** 3C cleavage site    **blue:** PIK3CB    **boxed:** E633K

PIK3CB wt <sup>1</sup> amino acid sequence							
1	<b>MCFSFIMPPA</b>	<b>MADILDIWAV</b>	<b>DSQIASDGS</b>	<b>PVDFLLPTGI</b>	<b>YIQLEVPREA</b>	<b>TISYIKQLMW</b>	60
61	<b>KQVHNYPMFN</b>	<b>LLMDIDSYMF</b>	<b>ACVNQTAVYE</b>	<b>ELEDETRRLC</b>	<b>DVRPFLPVLK</b>	<b>LVTRSCDPGE</b>	120
121	<b>KLDSKIGVLI</b>	<b>GKGLHEFDSL</b>	<b>KDPEVNEFR</b>	<b>KMRKFSEEKI</b>	<b>LSLVGLSWMD</b>	<b>WLKQTYPPEH</b>	180
181	<b>EPSIPENLED</b>	<b>KLYGGKLIVA</b>	<b>VHFENCQDVF</b>	<b>SFQVSPNMNP</b>	<b>IKVNELAIQK</b>	<b>RLTIHGKEDE</b>	240
241	<b>VSPYDYVLQV</b>	<b>SGRVEYVFGD</b>	<b>HPLIQFQYIR</b>	<b>NCVMNRALPH</b>	<b>FILVECKKIK</b>	<b>KMYEQEMIAI</b>	300
301	<b>EAAINRNSN</b>	<b>LPLPLPPKKT</b>	<b>RIISHVWENN</b>	<b>NPFQIVLVKG</b>	<b>NKLNTEETVK</b>	<b>VHVRAGLFHG</b>	360
361	<b>TELLCKTIVS</b>	<b>SEVSGKNDHI</b>	<b>WNEPLEFDIN</b>	<b>ICDLPRMARL</b>	<b>CFAVYAVLDK</b>	<b>VKTKKSTKTI</b>	420
421	<b>NPSKYQTIRK</b>	<b>AGKVHYPVAW</b>	<b>VNTMVFDFKG</b>	<b>QLRTGDIILH</b>	<b>SWSSFDPDELE</b>	<b>EMLNPMGTVO</b>	480
481	<b>TNPYTENATA</b>	<b>LHVKFPENKK</b>	<b>QPYYPFFDK</b>	<b>IEKAAEIAS</b>	<b>SDSANVSSRG</b>	<b>GKKFLPVLKE</b>	540
541	<b>ILDRDPLSQL</b>	<b>CENEMDLIWT</b>	<b>LRQDCREIFP</b>	<b>QSLPKLLLSI</b>	<b>KWNKLEDVAQ</b>	<b>LQALLQIWPK</b>	600
601	<b>LPPREALLEL</b>	<b>DFNYPDQYVR</b>	<b>EYAVGCLRQM</b>	<b>SDEELSQYLL</b>	<b>QLVQVLKYEP</b>	<b>FLDCALSRFL</b>	660
661	<b>LERALGNRRI</b>	<b>GQFLEWHLRS</b>	<b>EVHIPAVSQA</b>	<b>FGVILEAYCR</b>	<b>GSVGHMKVLS</b>	<b>KQVEALNKLK</b>	720
721	<b>TLNSLIKLN</b>	<b>VKLNRAKGKE</b>	<b>AMHTCLKQSA</b>	<b>YREALSDLQS</b>	<b>PLNPCVILSE</b>	<b>LYVEACKYMD</b>	780
781	<b>SKMKPLWLVA</b>	<b>NNKVFGEDEV</b>	<b>GVIKNGDSDL</b>	<b>RQDMLTLQML</b>	<b>RLMDLLWKEA</b>	<b>GLDLRMLPYG</b>	840
841	<b>CLATGDRSGL</b>	<b>IEVVSTSETI</b>	<b>ADIQLNSSNV</b>	<b>AAAAAFNKDA</b>	<b>LLNWLKEYNS</b>	<b>GDDLDRATIEE</b>	900
901	<b>FTLSCAGYCV</b>	<b>ASYVLGIGDR</b>	<b>HSDNIMVKKT</b>	<b>GQLFHIDFGH</b>	<b>ILGNFKSKFG</b>	<b>IKRERVPFIL</b>	960
961	<b>TYDFIHVIQQ</b>	<b>GKTGNTTEKFG</b>	<b>RFRQCCEADAY</b>	<b>LILRRHGNLF</b>	<b>ITLTFALMLTA</b>	<b>GLPELTSVKD</b>	1020
1021	<b>IQYKLDLAL</b>	<b>GKSEEEALKQ</b>	<b>FKQKFDALR</b>	<b>ESWTTKVNWM</b>	<b>AHTVRKDYRS</b>		1080

**blue:** PIK3CB sequence expressed in fusion protein    **Red:** variant in fusion protein

<sup>1</sup>NCBI/Protein accession number NP\_006210.1

Recombinant Proteins

PIK3R1 Recombinant Fusion Protein Amino Acid Sequence

1	MEEQKLISEE	DLPMVMSAEG	YQYRALYDYK	KEREEDIDLH	LGDILTVNKG	SLVALGFSDG	60
61	QEARPEEIGW	LNGYNETTGE	RGDFPGTYVE	YIGRKKISPP	TPKPRPPRPL	PVAPGSSKTE	120
121	ADVEQQALTL	PDLAEQFAPP	DIAPPLLIKL	VEAIEKKGLE	CSTLYRTQSS	SNLAE LRQLL	180
181	DCDTPSVDLE	MIDVHVLADA	FKRYLLDLPN	PVIPAAVYSE	MISLAPEVQS	SEEIYQLLKK	240
241	LIRSPSIPHQ	YWLTLQYLLK	HFFKLSQTSS	KNLLNARVLS	EIFSPMLFRF	SAASSDNTEN	300
301	LIKVIEILIS	TEWNERQPAP	ALPPKPPKPT	TVANNGMNNN	MSLQDAEWYW	GDISREEVNE	360
361	KLRDTADGTF	LVRDASTKMH	GDYTLTLRKG	GNNKLIKIFH	RDGKYGFSDP	LTFSSVVELI	420
421	NHYRNESLAQ	YNPKLDVKLL	YPVSKYQQDQ	VVKEDNIEAV	GKKLHRYNTQ	FQEKSRDYDR	480
481	LYEEYTRTSQ	EIQMKRTAIE	AFNETIKIFE	EQCQTQERYS	KEYIEKFKRE	GNEKEIQRIM	540
541	HNYDKLKSRI	SEIIDSRRRL	EEDLKKQAAE	YREIDKRMNS	IKPDLIQLRK	TRDQYLMWLT	600
601	QKGVRRQKLN	EWLGNENTED	QYSLVEDDED	LPHHDEKTWN	VGSSNRNKA	NLLRGKRDGT	660
661	FLVRESSKQG	CYACSVVVDG	EVKHCVINKT	ATGYGFAEPY	NLYSSLKELV	LHYQHTSLVQ	720
721	HNSLNVTLA	YPVYAQQRR					780

Red: MYC-tag blue: PIK3R1 K: E451K variation

PIK3R1 wt<sup>2</sup> amino acid sequence

1	MSAEGYQYRA	LYDYKKEREE	DIDLHLGDIL	TVNKGSLVAL	GFSDGQEARP	EEIGWLNQYN	60
61	ETTGERGDFP	GTVEYIIGRK	KISPPTPKPR	PPRPLPVAPG	SSKTEADVEQ	QALTLPLDLAE	120
121	QFAPPDIAPP	LLIKLVEAIE	KKGLECSTLY	RTQSSSNLAE	LRQLLDCDTP	SVDLEMIDVH	180
181	VLADAFKRYL	LDLNPVIPA	AVYSEMISLA	PEVQSSEEI	QLLKKLIRSP	SIPHQYWLTL	240
241	QYLLKHFFKL	SQTSSKNLLN	ARVLSEIFSP	MLFRFSAASS	DNTENLIKVI	EILISTEWNE	300
301	RQPAPALPPK	PPKPTTVANN	GMNNNMSLQD	AEWYWGDISR	EEVNEKLRDT	ADGTFLVRDA	360
361	STKMHGDYTL	TLRKGNNKL	IKIFHRDGKY	GFSDPLTFSS	VVELINHYRN	ESLAQYNPKL	420
421	DVKLLYPVSK	YQQDQVVKED	NIEAVGKKLH	EYNTQFQEK	REYDRLYEEY	TRTSQEIOMK	480
481	RTAIEAFNET	IKIFEEQCQT	QERYSKEYIE	KFKREGNEKE	IQRIMHNYDK	LKSRISEIID	540
541	SRRLEEDLK	KQAAEYREID	KRMNSIKPDL	IQLRKTRDQY	LMWLTQKGV	QKKLNEWLGN	600
601	ENTEDQYSLV	EDDEDLPHHD	EKTWNVGSSN	RNKAENLLRG	KRDGTFLVRE	SSKQGCYACS	660
661	VVDGGEVKHC	VINKTATGYG	FAEPYNLYSS	LKELVLHYQH	TSLVQHNSL	NVTLAYPVYA	720
721	QQRR						780

blue: PIK3R1 sequence expressed in fusion protein Red: variant in fusion protein

<sup>2</sup>NCBI/Protein accession number NP\_852664.1  
E451K: SNP variation see NCBI/dbSNP:rs17852841